



Glowing products for science

Isotype Control, Monoclonal Mouse IgG1 Kappa (IGG1/1331)

Non-reactive antibody to use as a negative control for mouse IgG1 primary antibodies. Available with CF® Dyes and other labels.



Product Description

This is a non-reactive mouse monoclonal IgG1 kappa antibody labeled with one of our bright and photostable CF® Dyes and other labels. This antibody may serve as a control for a mouse monoclonal IgG1 antibody under identical experimental conditions. Biotium also carries mouse monoclonal isotype control antibodies for [IgG2a Kappa \(MG2a-53\)](#) and [IgG2b Kappa \(IGG1/1333\)](#).

- Isotype control for antibody-based experiments
- Available with bright and photostable CF® Dyes
- APC, R-PE, RPE-Astral™616, RPE-Astral™775, APC-Astral™813, and biotin conjugates also available
- Suitable for western, flow cytometry, ELISA, immunofluorescence, and immunohistology

CF® Dyes offer exceptional brightness and photostability; learn more about [CF® Dyes](#).

Call us : [800-304-5357](tel:800-304-5357)

Product attributes

Antibody number	#1331
Antibody reactivity (target)	None
Antibody type	Isotype Control
Host species	Mouse
Clonality	Monoclonal
Clone	IGG1/1331
Isotype	IgG1, kappa
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Regulatory status	For research use only (RUO)
Antibody application notes	For coating for ELISA, order Ab without BSA, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry
Antibody/conjugate formulation	Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.
Species reactivity	None (isotype control)

Isotype Control, Monoclonal Mouse IgG1 Kappa (IGG1/1331)

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF@405S 1 mL (100 ug)	411/431 nm BNC041331-1ML	250 uL (25 ug)	BNC041331-250	CF@405S Features
CF@405M 1 mL (100 ug)	416/452 nm BNC051331-1ML	250 uL (25 ug)	BNC051331-250	CF@405M Features
CF@405L 1 mL (100 ug)	413/547 nm BNC061331-1ML	250 uL (25 ug)	BNC061331-250	CF@405L Features
CF@488A 1 mL (100 ug)	562/584 nm BNC881331-1ML	250 uL (25 ug)	BNC881331-250	CF@488A Features
CF@503R 1 mL (100 ug)	503/532 nm BNC501331-1ML	250 uL (25 ug)	BNC501331-250	CF@503R Features
CF@514 1 mL (100 ug)	516/549 nm BNC511331-1ML	250 uL (25 ug)	BNC511331-250	CF@514 Features
CF@568 1 mL (100 ug)	562/584 nm BNC681331-1ML	250 uL (25 ug)	BNC681331-250	CF@568 Features
CF@575 1 mL (100 ug)	575/601 nm BNC571331-1ML	250 uL (25 ug)	BNC571331-250	
CF@583R 1 mL (100 ug)	585/609 nm BNC581331-1ML	250 uL (25 ug)	BNC581331-250	CF@583R Features
CF@594 1 mL (100 ug)	593/615 nm BNC941331-1ML	250 uL (25 ug)	BNC941331-250	CF@594 Features
RPE-Astral™616 1 mL (100 ug)	496, 546, 566/617 nm BNC161331-1ML	250 uL (25 ug)	BNC161331-250	RPE-Astral™616 Features
CF@640R 1 mL (100 ug)	642/663 nm BNC401331-1ML	250 uL (25 ug)	BNC401331-250	CF@640R Features
CF@647 1 mL (100 ug)	652/668 nm BNC471331-1ML	250 uL (25 ug)	BNC471331-250	CF@647 Features
CF@647Plus 1 mL (100 ug)	652/668 nm BNC641331-1ML	250 uL (25 ug)	BNC641331-250	CF@647 Features
CF@660C 1 mL (100 ug)	667/685 nm BNC601331-1ML	250 uL (25 ug)	BNC601331-250	CF@660C Features
CF@700 1 mL (100 ug)	696/721 nm BNC001331-1ML	250 uL (25 ug)	BNC001331-250	CF@700 Features
CF@820 1 mL (100 ug)	822/835 nm BNC201331-1ML	250 uL (25 ug)	BNC201331-250	CF@820 Features
CF@870 1 mL (100 ug)	876/896 nm BNC701331-1ML	250 uL (25 ug)	BNC701331-250	CF@870 Features
RPE-Astral™775 1 mL (100 ug)	496, 546, 565/774 nm BNC751331-1ML	250 uL (25 ug)	BNC751331-250	RPE-Astral™775 Features
APC-Astral™813 1 mL (100 ug)	633, 638/813 nm BNC131331-1ML	250 uL (25 ug)	BNC131331-250	APC-Astral™813 Features
R-PE 1 mL (100 ug)	496, 546, 565/576 nm BNCR1331-1ML	250 uL (25 ug)	BNCR1331-250	
PerCP 250 uL (25 ug)	482/675 nm BNCP1331-250	1 mL (100 ug)	BNCP1331-1ML	
APC 1 mL (100 ug)	651/660 nm BNCA1331-1ML	250 uL (25 ug)	BNCA1331-250	
Biotin 1 mL (100 ug)	N/A BNCB1331-1ML	250 uL (25 ug)	BNCB1331-250	
Purified	N/A	250 uL (50 ug)	BNUB1331-250	
Purified, BSA-free	N/A	50 uL (50 ug)	BNUM1331-50	

This datasheet was generated on June 4, 2026 at 11:39:33 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/isotype-control-monoclonal-mouse-igg1-kappa-igg1-1331/>